

Prod. no. Cyp-01

Product Description
MOUSE ANTI-COD CYP1A
Monoclonal antibody, NP-7

biosense
LABORATORIES

www.biosense.com

Description

This product consists of affinity-purified mouse monoclonal antibodies against cytochrome P450 1A (CYP1A) from Atlantic cod. The antibody is dissolved in 0.1 M Na-phosphate, 0.15 M NaCl, 1% BSA, 0.1% Na-azide, pH 7.4.
Subclass/Isotype: IgG₁-kappa.

Immunogen source

Atlantic cod (*Gadus morhua* L.) cytochrome P450 1A, purified from liver microsomes of β -naphthoflavone treated cod.

Application

The monoclonal antibody Cyp-01 is well suited for a variety of applications including immunoblotting, ELISA and immunohistochemistry. Since assay conditions vary, the optimum dilution should be determined for each particular application.

Normal dilution range:

1:500 - 1:1 000 (Cod)

1:50 - 1:100 (Other species)

Storage

In the dark at 4°C.

Specificity

The monoclonal antibody Cyp-01 cross-reacts with a petroleum/PAH/PCB/dioxin-inducible protein ($m_r = 55,000-59,000$) in liver samples from a wide variety of different fish species including cod (*Gadiformes*), carp (*Cypriniformes*), salmon (*Salmoniformes*) and some flatfish (*Pleuronectiformes*) (see examples below). The degree of cross-reactivity differs between species and with the different methods used.

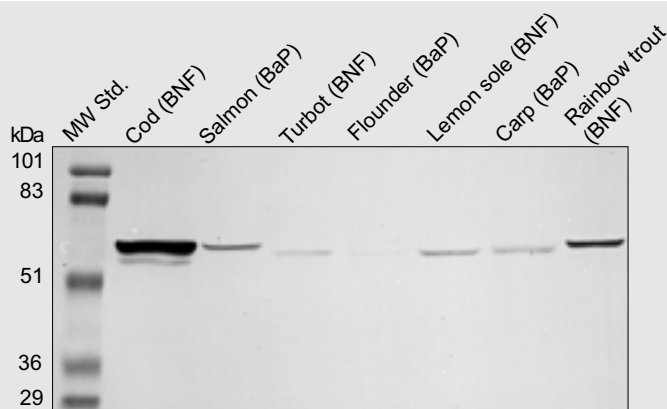
Western blot

Samples: Liver microsomes (10 μ g protein) from fish induced with β -naphthoflavone (BNF) or benzo(a)pyrene (BaP).

Primary antibody: Cyp-01 diluted 1:100

Secondary antibody: Peroxidase labelled Goat anti-mouse IgG, diluted 1:3000

Development: Peroxidase color development reagent (4-chloro-1-naphthol/ H_2O_2)



ELISA

Coating: Liver microsomes (1-10 μ g protein per well) from fish induced with β -naphthoflavone (BNF) or benzo(a)pyrene (BaP).

Primary antibody: Cyp-01

Secondary antibody: Peroxidase labelled Goat anti-mouse IgG, diluted 1:3000

Development: Peroxidase substrate (OPD/ H_2O_2)

